



MANAGEMENT SYSTEM CERTIFICATE

Certificate no.:
79976-2010-AQ-USA-ANAB

Initial certification date:
09 January, 2014

Valid:
20 November, 2024 – 11 December, 2025

This is to certify that the management system of
Tecan Precision Machining MH, Inc
885 Jarvis Drive, Morgan Hill, CA, 95037, USA
and the sites as mentioned in the appendix accompanying this certificate

has been found to conform to the Quality Management System standard:
ISO 9001:2015

This certificate is valid for the following scope:
The Provision of Machine Shop and Assembly Services

Place and date:
Katy, TX, 06 February, 2025

For the issuing office:
DNV - Business Assurance
1400 Ravello Drive, Katy, TX, 77449-5164, USA



Sherif Mekkawy
Management Representative



Certificate no.: 79976-2010-AQ-USA-ANAB
Place and date: Katy, TX, 06 February, 2025

Appendix to Certificate

Tecan Precision Machining MH, Inc

Locations included in the certification are as follows:

Site Name	Site Address	Site Scope
Tecan Precision Machining MH, Inc	885 Jarvis Drive, Morgan Hill, CA, 95037, USA	Management, Quality Assurance, Human Resources, Training, Competence, Production, Machine Shop, Inspection, Receiving, Stock, and Shipping
Tecan Precision Machining MH, Inc	875 Jarvis Drive, Morgan Hill, CA, 95037, USA	Assembly, Material Warehouse, Sales, Supply Chain Management including Outsourced Calibration

